

---

# LCK

Product code: 08-170

Full-length human LCK [1-509(end) amino acids of accession number NP\_005347.3] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 14 / 10

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 1.5

IC<sub>50</sub> at 1 mM ATP (nM) : 14

---